

---

# STREP A ISOLATION AGAR

---

## INTENDED USE

Remel Strep A Isolation Agar is a solid medium recommended for use in qualitative procedures for the primary isolation of beta-hemolytic group A streptococci from throat cultures and upper respiratory specimens.

## SUMMARY AND EXPLANATION

Pharyngitis caused by group A beta-hemolytic streptococci (*Streptococcus pyogenes*) can lead to serious medical complications, such as rheumatic fever and glomerulonephritis.<sup>1</sup> Throat culture is the primary means by which infection with *S. pyogenes* is identified. Several investigators have demonstrated improved isolation rates using sheep blood agar incorporated with selective agents to inhibit the commensal microbial flora commonly encountered in the oropharynx.<sup>2,3</sup> In 1977, Gunn et al. developed a selective medium consisting of sheep blood agar supplemented with sulfamethoxazole and trimethoprim for isolating groups A and B streptococci.<sup>4</sup> Using a modification of Gunn's formula, Welch et al. demonstrated superior recovery of *S. pyogenes* when compared with isolation rates found with nonselective sheep blood agar.<sup>5,6</sup>

## PRINCIPLE

Casein and soy peptones provide a nutritious source of organic nitrogen, amino acids, and peptides which are necessary for bacterial growth. Sodium chloride supplies essential electrolytes and maintains osmotic equilibrium. Sheep blood supplies nutrients necessary to support the growth of streptococci and allows detection of hemolytic reactions. A unique combination of selective ingredients is added to the medium to suppress commensal microbial flora and provide enhanced recovery of *S. pyogenes*.

## REAGENTS (CLASSICAL FORMULA)\*

Casein Peptone.....	15.0 g	Selective Agents.....	40.2 mg
Soy Peptone.....	5.0 g	Sheep Blood.....	5 %
Sodium Chloride.....	5.0 g	Agar.....	15.0 g
		Demineralized Water.....	1000.0 ml

pH 7.3 ± 0.2 @ 25°C

\*Adjusted as required to meet performance standards.

## PROCEDURE

1. Inoculate and streak the specimen as soon as possible after it is received in the laboratory. Selective and nonselective media should be inoculated to ensure recovery of microorganisms that may be inhibited on selective agar.
2. Inoculate Strep A Isolation Agar by rolling the swab over a small area of the agar surface. Use a sterile inoculating loop to streak the plate for isolation. Stab the agar several times with the loop in the area of heaviest inoculation. Anaerobiosis below the surface of the agar permits maximum expression of beta hemolysis by subsurface colonies.
3. Incubate plate in 5-10% CO<sub>2</sub> at 33-37°C for 24-48 hours.
4. Examine plate for typical colony morphology and beta hemolysis. On Strep A Isolation Agar, colonies of *S. pyogenes* are translucent or opaque, white to gray, and surrounded by a zone of beta hemolysis.

## QUALITY CONTROL

All lot numbers of Strep A Isolation Agar have been tested using the following quality control organisms and have been found to be acceptable. Testing of control organisms should be performed in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, patient results should not be reported.

### CONTROL

*Streptococcus pyogenes* ATCC® 19615  
*Escherichia coli* ATCC® 25922  
*Neisseria sicca* ATCC® 9913  
*Staphylococcus epidermidis* ATCC® 12228  
*Streptococcus pneumoniae* ATCC® 6305  
*Streptococcus sanguinis* ATCC® 10556

### INCUBATION

CO<sub>2</sub>, 18-24 h @ 33-37°C  
CO<sub>2</sub>, 18-24 h @ 33-37°C  
CO<sub>2</sub>, 18-24 h @ 33-37°C  
CO<sub>2</sub>, 18-24 h @ 33-37°C  
CO<sub>2</sub>, 18-24 h @ 33-37°C  
CO<sub>2</sub>, 18-24 h @ 33-37°C

### RESULTS

Growth, beta hemolysis  
Inhibition (partial to complete)  
Inhibition (partial to complete)  
Inhibition (partial to complete)  
Inhibition (partial to complete)  
Inhibition (partial to complete)

## PERFORMANCE CHARACTERISTICS

In a study of 600 throat swabs from patients suffering acute pharyngitis cultured on Strep A Isolation Agar and a standard, nonselective Blood Agar, the overall detection rates for all isolates of group A streptococci were 91% on Strep A Isolation Agar and 67% on nonselective Blood Agar. When both selective and nonselective agars were used in combination, the overall detection rates of group A streptococci increased to 93%.<sup>5</sup>

## LIMITATIONS

1. Colonies from selective media may not react with serogrouping reagents on direct testing. A subculture to a nonselective medium prior to serogrouping may be necessary.<sup>6</sup>
2. Organisms other than *S. pyogenes* (*Neisseria* spp., staphylococci, gram-negative bacilli) may grow on this medium but are generally inhibited. Additional biochemical and serological testing may be required for definitive identification. Consult appropriate references for further instructions.<sup>1</sup>

## BIBLIOGRAPHY

1. Murray, P.R., E.J. Baron, J.H. Jorgensen, M.L. Landry, and M.A. Pfaller. 2007. *Manual of Clinical Microbiology*. 9<sup>th</sup> ed. ASM Press, Washington, D.C.
2. Blanchette, L.P. and C. Lawrence. 1967. *Am. J. Clin. Pathol.* 48:441-443.
3. Vincent, W.F., W.E. Gibbons, and H.A. Gaafar. 1971. *Appl. Microbiol.* 22:942-943.
4. Gunn, B.A., D.K. Onashi, C.A. Gaydos, and E.S. Holt. 1977. *J. Clin. Microbiol.* 5:650-655.
5. Welch, D. 1989. Oklahoma Children's Memorial Hospital, Oklahoma City, OK. Data on file. Remel Inc., Lenexa, KS.
6. Welch, D.F., D. Hensel, D. Pickett, and S. Johnson. 1991. *Am. J. Clin. Pathol.* 95:587-590.

Refer to the front of Remel *Technical Manual of Microbiological Media* for **General Information** regarding precautions, product storage and deterioration, specimen collection, storage and transportation, materials required, quality control, and limitations.

ATCC® is a registered trademark of American Type Culture Collection.  
IFU 1856, Revised February 4, 2008

Printed in U.S.A.

# remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730    Technical Service: (800) 447-3641    Order Entry: (800) 447-3635

Local/International Phone: (913) 888-0939    International Fax: (913) 895-4128

Website: [www.remel.com](http://www.remel.com)    Email: [remel@remel.com](mailto:remel@remel.com)